



Instruction Manual

Catalog Number MU-0201

Version 01



Important Notice

Please, read the installation instruction carefully before installing the Smart Illuminator.

The Smart Illuminator is a LED illuminator intended for research use only. It must be operated only by qualified personnel aware of the potential health risks related to blue light irradiation and chemical and biological agents normally used in association with this device. Long-term exposure to blue light irradiation increases risk of retina diseases. The amber filter provides blue light protection for the end user but does not guarantee a complete protection.

Installation

Carefully unpack the Smart Illuminator as follow:

1. First remove the Smart Illuminator from its protective support.
2. Place it on a stable, horizontal surface.
3. Remove the plastic protection film from the transparent glass-viewing surface of the amber filter.

Stand alone installation

1. Keep at least 10 cm of space all around the Smart Illuminator to ensure good ventilation.
2. Use the supplied cable to connect the device to the power. The power font must be able to deliver at least 250 VA with a voltage between 100 and 240 Vac. The plug must have a ground connection. The adaptor supplies 24 V DC voltage.

General Precautions

- Plug the Smart Illuminator on an electric line with ground connection.
- The Smart Illuminator is equipped with thermal protection to prevent overheating.
- Do not pour liquids directly on the illuminator.
- Do not block the aeration slits.
- Switch off the instrument immediately after its use.
- Position the Smart Illuminator to prevent harm to nearby operators.
- The Smart Illuminator sample surface is a glass sheet for cutting gels. When using the illuminator with samples stained with specific dyes, decontaminate the illuminator surface with bleach. Denatured alcohol can be also used. Always wear disposable gloves.

Using the Smart Illuminator

1. Place gel/sample on the transparent glass support area.
 Please wear gloves to prevent skin contact with gel and staining agents.
2. Press the ON/OFF switch to ON to turn the LEDs on.
 Please observe the gel through the amber filter. Long-term exposure to blue light irradiation increases risk of retina diseases.
3. After viewing the sample or cutting out the bands, turn the illuminator off.
 Continuous or frequent LED lighting during electrophoresis of gels may decay the DNA fluorescence intensity.

General Appearance of the Smart Illuminator



Specifications

- Dimensions (mm): 210D x 210W x 30H
- Viewing surface (mm): 120 x 70
- Wavelength (nm): 470
- Amber filter: amber filter shield with metal frame
- LED arrangements: matrix for two-side illumination
- LED lifetime: 50,000 hours
- Power: 24V DC 1.25 A power adaptor
- Compatible with mini gel size (mm): 110 x 60 & 55 x 60
- Weight (Kg): 2.3

Front Panel of the Smart Illuminator



Rear Panel of the Smart Illuminator



Adaptor: 24V DC / 1.25 A

Power AC input: 100~240 V; 50/60 Hz power source

Warranty

The Smart Illuminator is warranted against defects in materials and workmanship for 2 year. If any defects occur in the instrument or accessories during this warranty period, Eurogentec will repair or replace the defective parts at its discretion without any charge. The following defects, however, are specifically excluded:

1. Defects caused by improper operation.
2. Repair or modification done by anyone other than Eurogentec or an authorized agent.
3. Damage caused by substituting alternative parts.
4. Use of fittings or spare parts supplied by anyone other than Eurogentec.
5. Damage caused by accident or misuse.
6. Damage caused by disaster.
7. Corrosion caused by improper solvent or sample.

For any inquiry or request for repair service, please contact our technical support at:
customer.care@eurogentec.com

Regulatory Notice

The Smart Illuminator is designed and certified to meet safety standards and EMC regulations. Certified products are safe to use when operated in accordance with the instruction manual. This instrument should not be modified or altered in any way. Alteration of this instrument will:

1. Void the manufacturer's warranty
2. Void the safety and EMC certification
3. Create a potential safety hazard

Eurogentec is not responsible for any injury or damage caused by the use of this instrument for purposes other than those for which it is intended, or by modifications of the instrument not performed by Eurogentec or an authorized agent.

CONTACT

EU : Eurogentec Headquarters, Europe
email : info@eurogentec.com
Toll Free Number : 00 800 666 00 123

US : AnaSpec Inc., North America
email : service@anaspec.com
Toll Free Number : +1 800 452 5530

Eurogentec is part of Kaneka Corporation

